

Mastoparan

Mast cell degranulating peptide

Cell permeable, amphiphilic, mast cell degranulating tetradecapeptide. Stimulates glycogenolysis and phospholipase D₂ (PLD₂) activity. Alters G-proteins in brain membranes. Directly activates G proteins by a mechanism analogous to that of G protein-coupled receptors. Acts preferentially on G_i and G_o rather than G_s. Stimulates exocytosis and phosphoinositide breakdown in a variety of cells. Inhibits calmodulin, and activates PLA₂. Acts as a mitogen in Swiss 3T3 cells and stimulates pertussis toxin-sensitive arachidonate release without phosphoinositide breakdown.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-G410-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	72093-21-1
Formula	$C_{70}H_{131}N_{19}O_{15}$
Formulation	Lyophilized.
Identity	Determined by MS.
MW	1478.9
Purity	≥95% (HPLC)
Sequence	H-Ile-Asn-Leu-Lys-Ala-Leu-Ala-Ala-Leu-Ala-Lys-Lys-Ile-Leu-NH ₂
Solubility	Soluble in water (1 mg/ml) or 100% ethanol.
Source	Synthetic. Originally isolated from wasp venom (<i>Vespula lewisii</i>).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)